Complement component C1s (human) monoclonal antibody (9)

C1s proenzyme is a 79.8kDa glycoprotein produced primarily by hepatocytes but also by other cells. The proenzyme is produced as a single-chain polypeptide of 688 amino acids. Upon activation, the peptide is cleaved between R437 and I438, producing an activated enzyme consisting of two disulfide-linked chains, A chain (52.2kDa) and B chain (27.7kDa). The calcium-dependent catalytic subunit of the C1 complex (C1q, C1r, C1s) is comprised of two C1s and two C1r molecules, and this serine protease tetramer is responsible for the initial steps in the cleavage of C4 and C2 in the classical complement pathway.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-ABS-002-09-1	1mg
BPD-ABS-002-09-04	400µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name C1 esterase, Complement C1s subcomponent, C1s

proenzyme

Application ELISA, IHC

Clone 9

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

GenBank ID 716

Host Mouse

ImmunogenHuman plasma complement component C1s Proenzyme.

Isotype IgG1ĸ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

Specificity Specific for C1s proenzyme and activated free C1s.

UniProt ID P09871

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com

Last modified: May 29, 2024